

Anti-human IFN- α

Mouse Monoclonal Antibody

Cat. No.	Form	Volume	Quantity
BMS160	Pure	100 μ l	100 μ g

Clone	EBI-1
Subclass	mouse IgG1Kappa
Purity	BMS160 was purified by ammonium sulfate precipitation and DEAE sepharose chromatography to >90% purity as assessed by SDS-PAGE.
Presentation	PBS
Specificity	BMS160 specifically reacts with human recombinant and natural interferon - α
Applications	<p>ELISA: BMS160 can be used as capture antibody in ELISA.</p> <p>Functional studies: BMS160 inhibits the antiviral activity of interferon -α. Optimal dilutions should be determined by the individual laboratory for each application.</p>
Cross-reactivity	BMS160 does not crossreact with human IFN- ω , IFN- α B, IFN- α F.
Antigen	Recombinant human interferon- α 2 $^{\circ}$
Storage	2-8 $^{\circ}$ C until expiry as indicated on label. After opening aliquote contents and freeze at -20 $^{\circ}$ C.
Shipping conditions	2-8 $^{\circ}$ C

Limited product warranty

Life Technologies Corporation and/ or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at thermofisher.com/support.

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

Corporate entity: Life Technologies | Carlsbad, CA 92008 USA | Toll Free in USA 1.800.955.6288

©2017 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. All other trademarks are properties of their respective owners.

For Research Use Only. Not for use in diagnostic procedures.